

---

---

# MAP3K1

Product code: 07-103

Human MAP3K1, catalytic domain [1193-1512(end) amino acids of accession number NP\_005912.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : MAP2K1  
Cascade Assay \*  
ATP ( $\mu$ M) : 1000  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at 1 mM ATP (nM) : 160

\* MAP2K1/Erk2/Modified Erktide

---

---